

Rabbit Anti-Alexa Fluor 350 Antibody, iFluor® 555-Labeled

Catalog Number: 46628

Unit Size: 100 ug

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Liquid
Molecular Weight	N/A
Soluble In	PBS

Spectral Properties

Excitation Wavelength	557 nm
Emission Wavelength	570 nm

Applications

Rabbit Anti-Alexa Fluor 350 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 350-conjugated molecules and is directly labeled with iFluor® 555, a bright orange fluorophore with excitation/emission at around 555/565 nm. This direct labeling format enables one-step fluorescence detection without the need for a secondary antibody.

The antibody may cross-react with related blue fluorophores such as iFluor® 350, AMCA, and Cascade Blue®. Its orange emission is spectrally distinct from Alexa Fluor® 350, making it ideal for multicolor imaging, flow cytometry, and other fluorescence-based assays.

Alexa Fluor® and Cascade Blue® are registered trademarks of Thermo Fisher Scientific.